

Product datasheet

Dihydroxyacetone Phosphate (DHAP) Assay Kit (Fluorometric) ab197003

[3 References](#) [2 Images](#)

Overview

Product name	Dihydroxyacetone Phosphate (DHAP) Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Serum, Plasma, Other biological fluids, Tissue, Adherent cells, Suspension cells
Assay type	Quantitative
Sensitivity	> 0.5 μ M
Species reactivity	Reacts with: Mammals, Other species
Product overview	Dihydroxyacetone Phosphate (DHAP) Assay Kit (Fluorometric) (ab197003) is suitable for measuring low levels of DHAP typically found in a variety of samples. In this kit, triose phosphate isomerase (TPI) converts dihydroxyacetone phosphate (DHAP) to glyceraldehyde-3-phosphate (GAP), which subsequently undergoes a series of reaction and reduces the probe to generate fluorescence. The fluorescence intensity generated is directly proportional to the amount of dihydroxyacetone Phosphate.

Detection limit: 0.5 μ M DHAP.

Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K673 PicoProbe™ Dihydroxyacetone Phosphate (DHAP) Fluorometric Assay Kit. K673-100 is the same size as the 100 test size of ab197003.</p> <p>Dihydroxyacetone Phosphate (DHAP) is an important intermediate in both lipid biosynthesis and glycolysis. In glycolysis, fructose-1,6- diphosphate is converted to dihydroxyacetone phosphate (DHAP) and glyceraldehyde-3-phosphate (GAP) by aldolase. Both DHAP and GAP serve as the intracellular pool for triose phosphate. DHAP can be further converted into GAP by Triose Phosphate Isomerase (TPI).</p> <p>In humans, TPI deficiency is a rare autosomal disease. It causes hemolytic anemia, neurological diseases, and even death due to blockage of the glycolytic pathway and accumulation of DHAP in erythrocytes.</p>
--------------	--

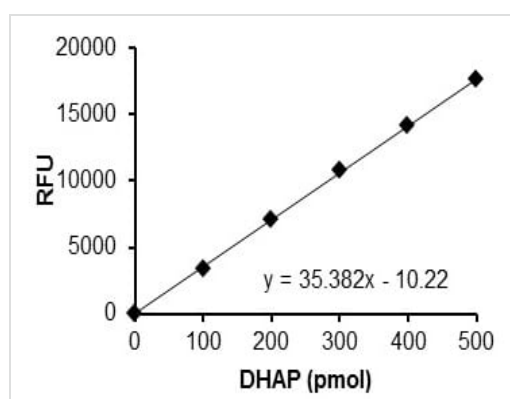
Platform	Microplate reader
-----------------	-------------------

Properties

Storage instructions	Store at -20°C. Please refer to protocols.
-----------------------------	--

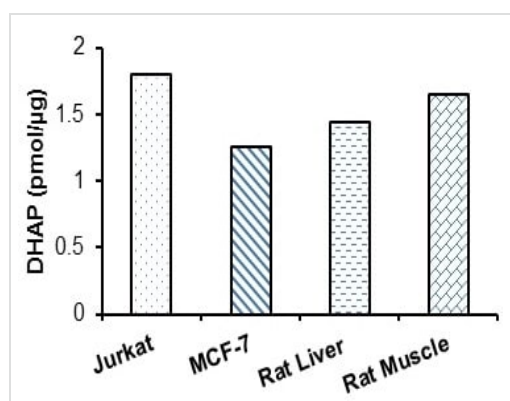
Components	100 tests
Assay Buffer II	1 x 25ml
Developer Solution X	1 vial
DHAP Enzyme Mix	1 vial
DHAP Standard	1 vial
PicoProbe I	1 x 400µl

Images



Typical DHAP Standard Curve obtained following assay protocol.

Functional Studies - Dihydroxyacetone Phosphate
(DHAP) Assay Kit (Fluorometric) (ab197003)



Measurement of DHAP level in a variety of samples: Jurkat (250 µg), and MCF-7 (150 µg) cell lysate, and in rat liver (50 µg) and rat muscle (150 µg) tissue lysate.

Functional Studies - Dihydroxyacetone Phosphate
(DHAP) Assay Kit (Fluorometric) (ab197003)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors